

# Benserazide . HCl

## Peripheral decarboxylase inhibitor

Peripheral decarboxylase inhibitor. Also inhibits tryptophan hydroxylase and tyrosine hydroxylase. Used with L-DOPA as a Parkinson's disease therapeutic.

Citations: 6

[View Online »](#)

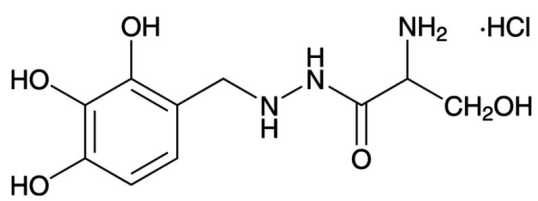
## Ordering Information

[Order Online »](#)

|                |    |
|----------------|----|
| BML-EI165-1000 | 1g |
|----------------|----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |   |
|-------------------|---|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Product is subject to oxidation. Make solutions fresh daily. Store, as supplied, at room temperature for up to 1 year. |
| Handling          | Protect from light.   |
| Long Term Storage | Ambient   |
| Shipping          | Ambient Temperature   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|               |   |
|---------------|---|
| Appearance    | Tan solid.  |
| CAS           | 14919-77-8  |
| Couple Target | Peripheral decarboxylase, Tryptophan hydroxylase, Tyrosine hydroxylase    |
| Couple Type   | Inhibitor   |
| Formula       | $C_{10}H_{15}N_3O_5 \cdot HCl$  |
| MW            | 293.7   |
| Purity        | ≥98%  |
| Solubility    | Soluble in water. Almost insoluble in 100% ethanol, insoluble in acetone. |



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)